

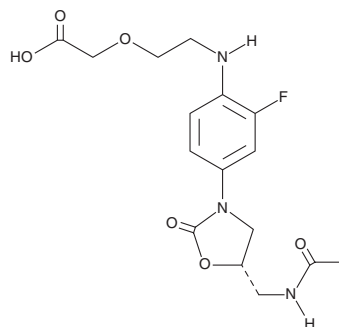
PRODUCT INFORMATION



PNU 142300

Item No. 26608

CAS Registry No.: 368891-69-4
Formal Name: 2-[2-[[4-[(5S)-5-[(acetylamino)methyl]-2-oxo-3-oxazolidinyl]-2-fluorophenyl]amino]ethoxy]-acetic acid
MF: C₁₆H₂₀FN₃O₆
FW: 369.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PNU 142300 is supplied as a solid. A stock solution may be made by dissolving the PNU 142300 in the solvent of choice, which should be purged with an inert gas. PNU 142300 is slightly soluble in the organic solvent DMSO. PNU 142300 is also slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

PNU 142300 is an inactive metabolite of the oxazolidinone antibiotic linezolid (Item No. 15012).¹

Reference

1. Slatter, J.G., Stalker, D.J., Feenstra, K.L., *et al.* Pharmacokinetics, metabolism, and excretion of linezolid following an oral dose of [(¹⁴C)]linezolid to healthy human subjects. *Drug. Metab. Dispos.* **29(8)**, 1136-1145 (2001).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/29/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM